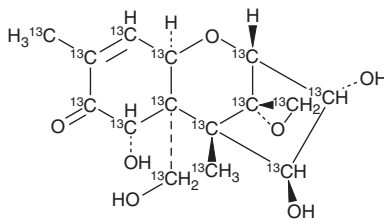


PRODUCT INFORMATION



$^{13}\text{C}_{15}$ -Nivalenol
Item No. 31259

CAS Registry No.: 911392-40-0
Formal Name: (3 α ,4 β ,7 α)-12,13-epoxy-3,4,7,15-tetrahydroxy-trichothec-9-en-8-one- $^{13}\text{C}_{15}$
MF: [$^{13}\text{C}_{15}$]H₂₀O₇
FW: 327.2
Purity: $\geq 98\%$
Supplied as: A 25 $\mu\text{g}/\text{ml}$ solution in acetonitrile
Storage: -20°C
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

$^{13}\text{C}_{15}$ -Nivalenol is intended for use as an internal standard for the quantification of nivalenol (Item No. 11438) by GC- or LC-MS. Nivalenol is a trichothecene mycotoxin that has been found in *Fusarium*.¹ It is lethal to mice ($\text{LD}_{50} = 6.9 \text{ mg}/\text{kg}$).² Nivalenol (5, 10, and 15 mg/kg) also induces thymic, splenic, and Peyer's patch cell apoptosis in mice.³

References

1. Yang, Z., Concannon, J., Ng, K.S., *et al.* Tetrandrine identified in a small molecule screen to activate mesenchymal stem cells for enhanced immunomodulation. *Sci. Rep.* **6**, 30263 (2016).
2. Yoshizawa, T. and Morooka, N. Studies on the toxic substances in the infected cereals (part 3): Acute toxicities of new trichothecene mycotoxins: Deoxynivalenol and its monoacetate. *J. Food Hyg.* **15**(4), 261-269 (1974).
3. Poapolathep, A., Ohtsuka, R., Kiatipattanasakul, W., *et al.* Nivalenol-induced apoptosis in thymus, spleen and Peyer's patches of mice. *Exp. Toxicol. Pathol.* **53**(6), 441-446 (2002).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/09/2026

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM